

PRODUCT INFORMATION



1-Myristoyl-2-Linoleoyl-3-Palmitoyl-*rac*-glycerol

Item No. 28560

CAS Registry No.: 105595-41-3

Formal Name: 9Z,12Z-octadecadienoic acid,
1-[[[(1-oxohexadecyl)oxy]methyl]-
2-[[[(1-oxotetradecyl)oxy]ethyl ester
Synonyms: 1-Myristin-2-Linolein-3-Palmitin,
TG(14:0/18:2/16:0)

MF: $C_{51}H_{94}O_6$

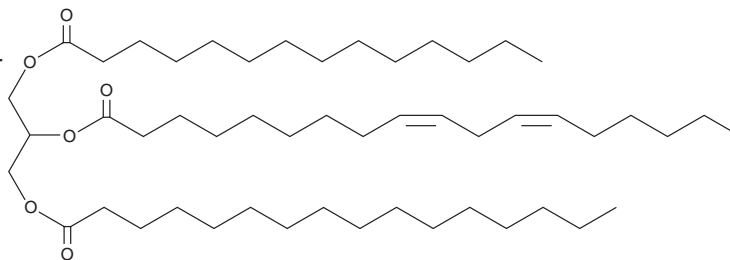
FW: 803.3

Purity: $\geq 95\%$

Supplied as: A solution in methyl acetate

Storage: -20°C

Stability: ≥ 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

1-Myristoyl-2-linoleoyl-3-palmitoyl-*rac*-glycerol is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. 1-Myristoyl-2-linoleoyl-3-palmitoyl-*rac*-glycerol is slightly soluble in chloroform and methanol.

Description

1-Myristoyl-2-linoleoyl-3-palmitoyl-*rac*-glycerol is a triacylglycerol that contains myristic acid (Item No. 13351), linoleic acid (Item Nos. 90150 | 90150.1 | 21909), and palmitic acid (Item No. 10006627) at the *sn*-1, *sn*-2, and *sn*-3 positions, respectively. It has been found in palm and vegetable oils.^{1,2}

References

1. Sassano, G.J. and Jeffrey, B.S.J. Gas chromatography of triacylglycerols in palm oil fractions with medium-polarity wide-bore columns. *J. Am. Oil Chem. Soc.* **70**(11), 1111-1114 (1993).
2. Ohshima, T., Yoon, H.-S., and Koizumi, C. Application of selective ion monitoring to the analysis of molecular species of vegetable oil triacylglycerols separated by open-tubular column GLC on a methylphenylsilicone phase at high temperature. *Lipids* **24**(6), 535-544 (1989).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/22/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM